

Custom oligonucleotides Additional Services	
Item Number	Additional Services for custom oligonucleotides Design requests
AD-DGDDP	Double Dye Probe design (including LNA®)
AD-DGDDP-SNP	Double Dye Probe design (including LNA®) for SNP Analysis
AD-DGDSP	Duplex Scorpions® Probes design
AD-DGDSP-SNP	Duplex Scorpions® Probes design for SNP Analysis
AD-DGMBN	Molecular Beacon design
AD-DGMBN-SNP	Molecular Beacon design for SNP Analysis
AD-DGPCR	PCR Primers (Forward + Reverse)
AD-DGSBE-SNP	Scorpions® Probe design for SNP Analysis
AD-DGSPE	Scorpions® Probe design
Additional Services for custom oligonucleotides Format options	
AD-FRANN-001	Annealed as Duplex
AD-FRMIX-001	Mixed Oligonucleotides
AD-FRSOL-001	In Solution (H ₂ O, TE...)
Additional Services for custom oligonucleotides Packaging options	
AD-PK096-CLUS	96-Well Plate Cluster Tubes
AD-PK096-DEEP	96-DeepWell Plate
AD-PK096-WELL	96-Well Plate
AD-PK384-GREIN	384-Well Plate
AD-PKALQ-001	Aliquoting in X tubes
Additional Services for custom oligonucleotides Quality Controls	
AD-QCCGE-001	Capillary Gel Electrophoresis (CGE)
AD-QCFLU-001	Fluorescence Analysis
AD-QCHPL-IEX	Ion-Exchange HPLC (IEX-HPLC)
AD-QCHPL-RP	Reverse Phase HPLC (RP-HPLC)
AD-QCMMS-001	MALDI-TOF MS
AD-QCMMS-002	MALDI-TOF MS Sequencing (Non-Modified Oligonucleotide)
AD-QCMMS-003	MALDI-TOF MS Sequencing (Modified Oligonucleotide)
AD-QCMMS-004	MALDI-TOF MS Sequencing (Double-Dye Probe)